

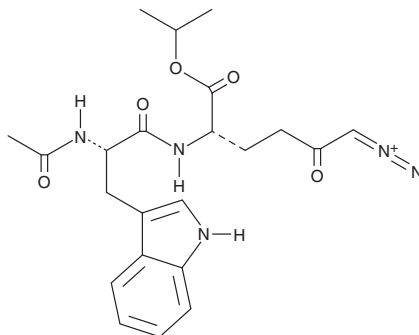
PRODUCT INFORMATION



Sirpiglenastat

Item No. 41293

CAS Registry No.: 2079939-05-0
Formal Name: N-acetyl-L-tryptophyl-6-diazo-5-oxo-L-norleucine, 1-methylethyl ester
Synonym: DRP-104
MF: C₂₂H₂₇N₅O₅
FW: 441.5
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Sirpiglenastat is supplied as a solid. A stock solution may be made by dissolving the sirpiglenastat in the solvent of choice, which should be purged with an inert gas. Sirpiglenastat is sparingly soluble (1-10 mg/ml) in DMSO.

Description

Sirpiglenastat is a prodrug form of the glutaminase inhibitor 6-diazo-5-oxo-L-nor-leucine (DON; Item No. 17580).¹ Sirpiglenastat decreases P493B lymphoma cell viability (EC₅₀ = 10 μM). *In vivo*, sirpiglenastat is metabolized to DON in tumor tissues, but metabolized to an inactive metabolite in gastrointestinal tissues, in EL4 murine T cell lymphoma, or MC-38 murine colon, EO771 murine breast, or 3LL Lewis lung carcinoma tumor-bearing CES1^{-/-} mice, a model designed to more closely mimic human metabolism. It reduces tumor size in EL4 tumor-bearing CES1^{-/-} mice when administered at doses of 0.1, 0.3, and 1 mg/kg.

Reference

1. Rais, R., Lemberg, K.M., Tenora, L., *et al.* Discovery of DRP-104, a tumor-targeted metabolic inhibitor prodrug. *Sci. Adv.* **8**(46), eabq5925 (2022).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 01/10/2025

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM